

Data Sheet

TIGIT, Fc Fusion (IgG1), Avi-Tag, Biotin (Mouse)

Mouse, Recombinant, Fc fusion protein,
 N-terminal Avi-Tag, Biotin-Labeled
Catalog #: 79269
Lot #: 170824 **Conc.:** 0.75 mg/ml

Formulated in: 8mM phosphate, pH7.4,
 110mM NaCl, 2.2mM KCl, and 20% glycerol.

Stability: At least 6 months at -80°C . Avoid
 freeze/thaw cycles. Protein may be diluted to
 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at
 -80°C .

References:

1. Yu, X., *et al.*, *Nat. Immunol.* 2009; **10(1)**:
 48-57.
2. Stanietsky, N., *et al.*, *Proc. Natl. Acad.
 Sci. (USA)*. 2009; **106(42)**: 17858-17863.

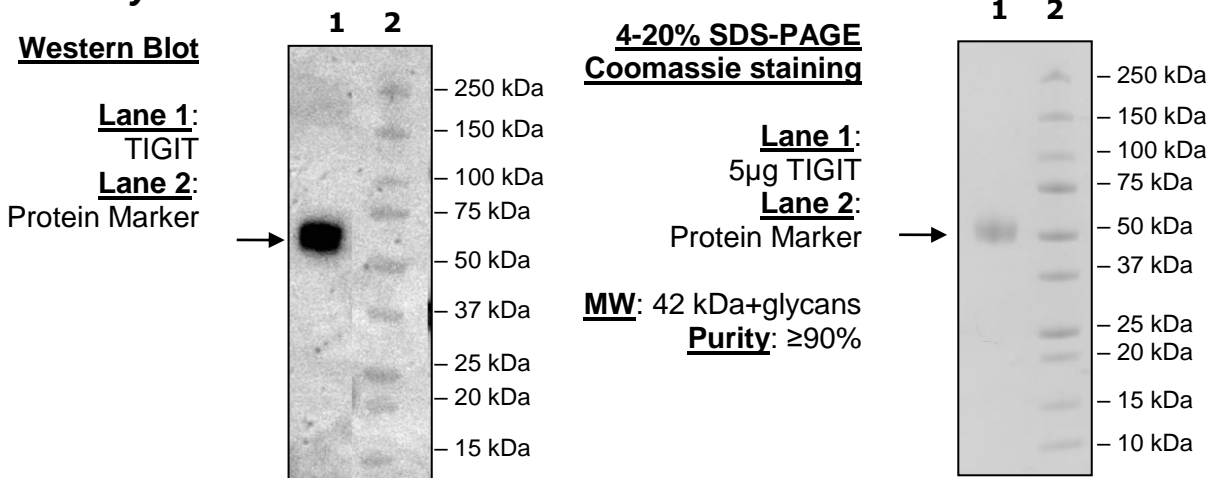
Description:

Mouse T-cell immunoreceptor with Ig and
 ITIM domains (TIGIT), also known as V-set
 and immunoglobulin domain-containing
 protein 9, VSIG9, V-set and transmembrane
 domain-containing protein 3, and VSTM3,
 GenBank Accession No. NM_001146325.1,
 a.a. 26-143 fused to Fc region of human
 IgG, expressed in a HEK293 cell expression
 system. MW = 42 kDa. This protein runs at
 a higher MW by SDS-PAGE due to
 glycosylation.

Application:

Useful as a positive control for western
 blotting.

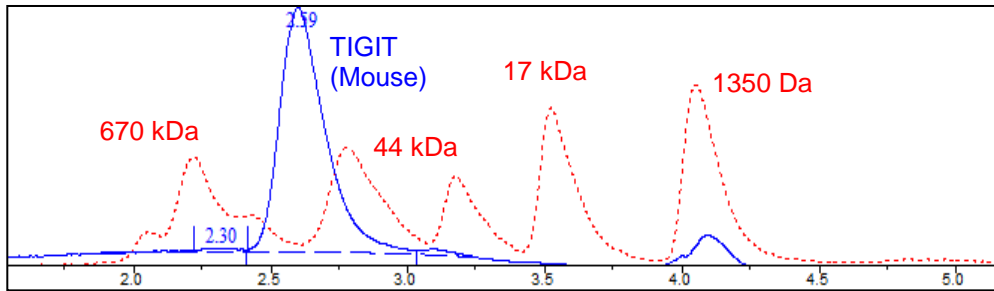
Quality Assurance



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com

Gel Filtration Curve



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com